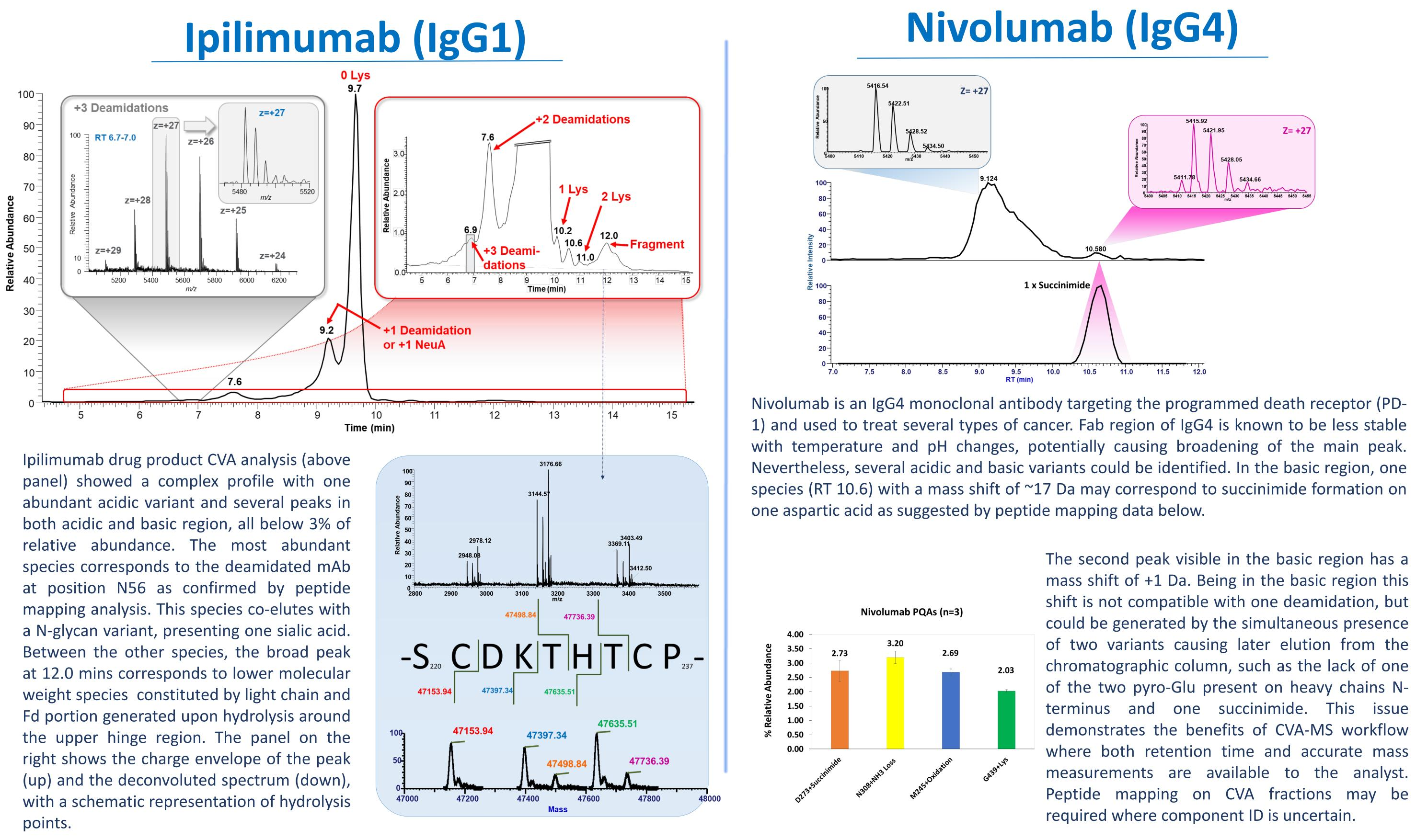
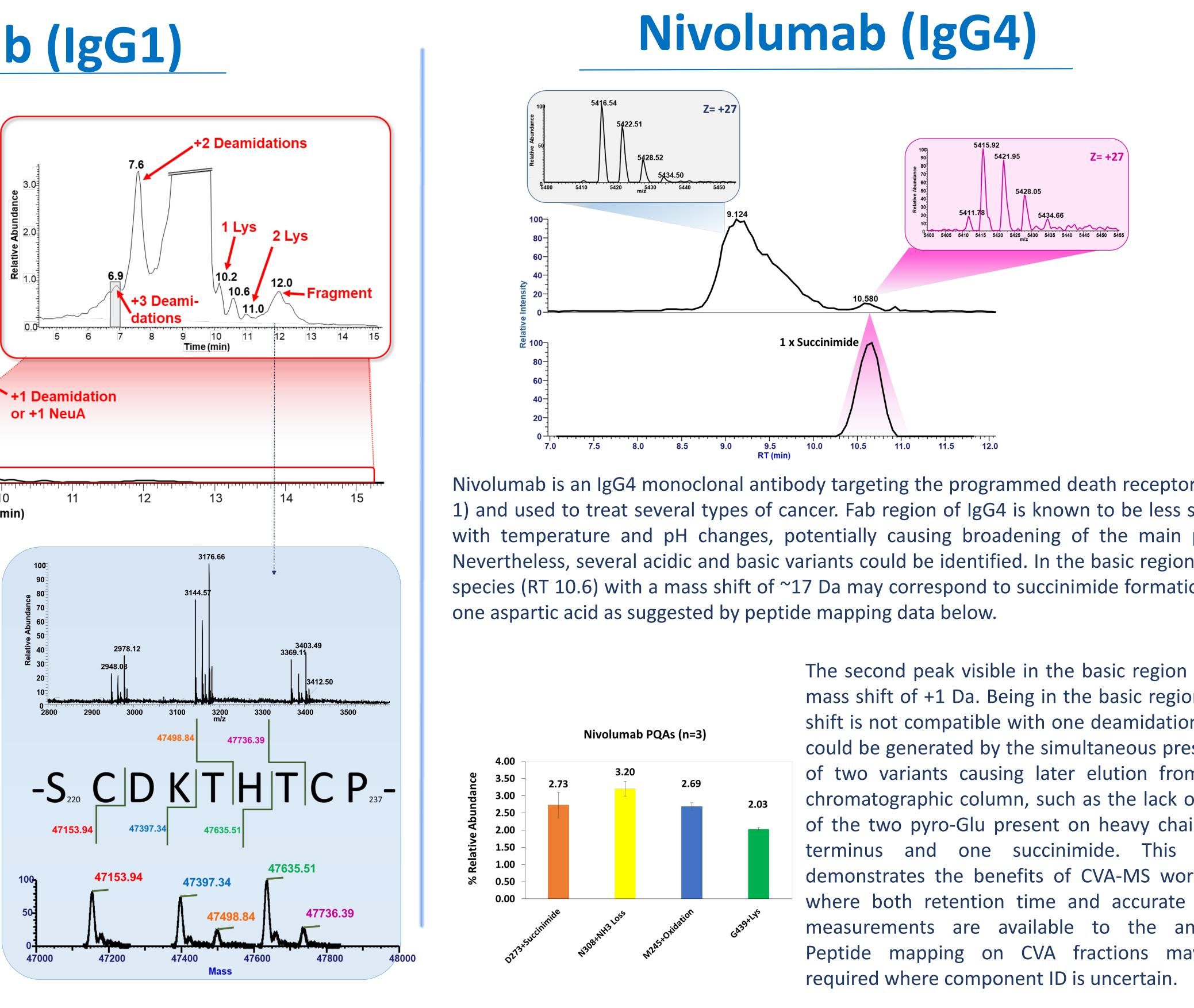
## Monoclonal antibody characterization through native Orbitrap mass spectrometry leading to improved sensitivity and elucidation of microheterogeneity

<u>Sara Carillo<sup>1</sup>; Silvia Millán Martín<sup>1</sup>; Florian Füssl<sup>1</sup>; Itzcoatl Gomez Aquino<sup>2</sup>; Ioscani Jimenez Del Val<sup>2</sup>; Jonathan Bones<sup>1, 2</sup></u> <sup>1</sup>National Institute for Bioprocessing Research and Training (NIBRT), Blackrock, Ireland; <sup>2</sup>University College of Dublin, Dublin, Ireland



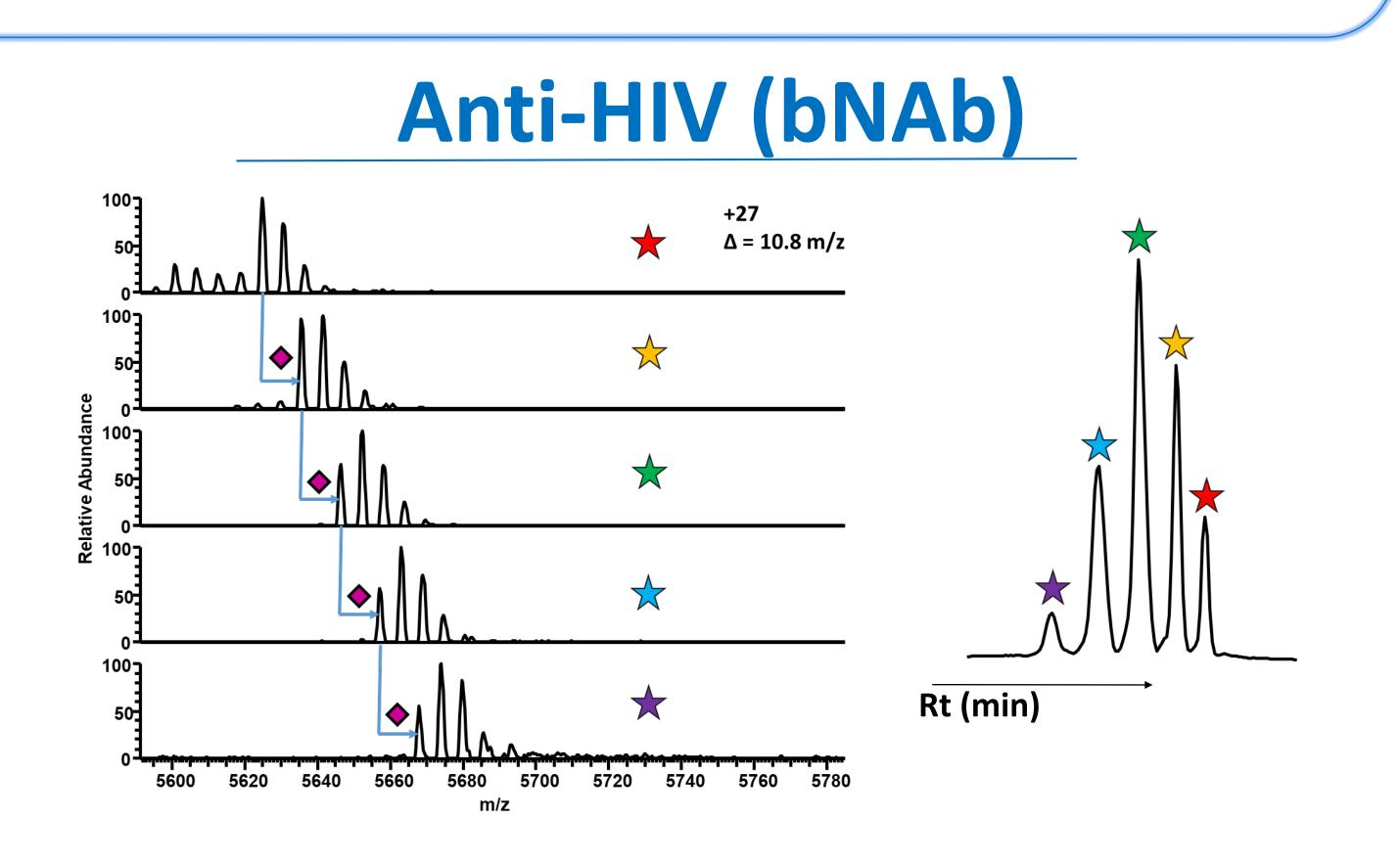
- window option using ReSpect<sup>™</sup> algorithm.





Conclusions: In the present study 3 mAbs were analysed using CVA analysis hyphenated with new generation Orbitrap mass spectrometry instrument. For all samples, Acknowledge Thermo Fisher Scientific for instrument access and support. excellent data quality and identification of low abundant variants within 25 ppm was achieved also on very low abundant proteoforms (< 1%). Charge variant analysis on the References: [1] Tassi, M. et al. Journal of Separation Science 41, 125-144, (2017). [2] Füssl, F. et al. Analytical Chemistry Orbitrap Exploris 240 MS enables in-depth characterization and confident identification of mAb microheterogeneity due to excellent sensitivity and spectral quality. 90 (7), 4669-4676, (2018). [3] Füssl, F. et al. MABS, 11 (1), 116-128, (2018).

Background: Mass spectrometry (MS) of intact proteins is increasingly used in biopharmaceutical analysis as it is rapid and provides significant structural insights without laborious sample preparation steps that may interfere with endogenous post-translational modifications (PTMs) present on the drug substance. Significant advances in the use of chromatography under native conditions and the introduction of volatile mobile phases [1]. Charge variant analysis (CVA) can be performed using ion exchange chromatography. Using volatile buffers with low salt concentrations it is possible to provide compatibility with MS, enabling detailed characterization of low abundant modifications that may be more difficult to obtain with size exclusion chromatography (SEC) [2,3]. Sensitivity and performance at higher resolution settings, may provide increased dynamic range and confidence in charge variant identification. Orthogonal techniques, such as peptide mapping and subunit analysis, may be employed to confirm the identity of some very low abundant species. **Methodology:** CVA analysis was performed using a Thermo Scientific<sup>™</sup> MAbPac<sup>™</sup> SCX-10 RS (2.1 x 50 mm) column. Linear gradient using A) 25 mM ammonium bicarbonate, 30 mM acetic acid (pH 5.3) and B) 10 mM ammonium hydroxide (pH 10.9). Separation was performed on a Vanquish<sup>™</sup> UHPLC (Thermo Scientific) hyphenated with an Thermo Scientific<sup>™</sup> Orbitrap Exploris<sup>™</sup> 240 mass spectrometer. Sample amount: 50 µg injected in triplicate. Data processing was performed using BioPharma<sup>™</sup> Finder software v. 4.0, Sliding



The third antibody analysed is a broadly neutralising antibody (bNAb) targeting HIV and produced in CHO cell line. A second N-glycosylation site is present on bNAb light chain, with sialylated non-fucosylated complex N-glycans, which influence CVA profile at intact level (panel above). Only performing CVA analysis on the Fab region obtained after IdeS digestion (lower panel) it was possible to observe fucose presence on the N-glycans present on the light chain.

